

Instruction Manual

Catalog Number	PK-AB577-3606
Description	Human HDAC-6 (1215 a.a. residues) possesses two separate putative catalytic domains, both of which are fully functional and contribute independently to the overall activity of HDAC-6. A very potent NES is present at the amino-terminus of HDAC-6, which was found to play an important role in regulating the shuttling of HDAC-6 protein between cytoplasm and nucleus. The shuttling process may be a critical regulatory mechanism of HDAC-6 function. HDAC-6 may participate in coordinating expression of a group of genes involved in the remodeling of chromatin during cell differentiation.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide mapping to the N-terminus of mouse HDAC-6 (ID BV-48).
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	The antibody detects ~134 kDa HDAC-6. The antibody does not cross-react with other HDAC proteins including HDAC-1, 2, 3, 4, 5, 7, and 8.
Formulation	100 µg (0.2 mg/ml) affinity purified rabbit anti-HDAC-6 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and freeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	The antibody can be used in Western blotting (0.5-4 µg/ml). Based on researcher's feedback, the antibody can also be used in Immunohistochemistry (20 µg/ml). However, the optimal conditions should be determined individually. Jurkat cell lysate (Cat.# PK-AB718-1205) can be used as a positive control. Blocking peptide (Cat.# PK-AB577-3606P) is available separately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.